

Human SAA ELISA Kit (SAA1) ab100635

[13 References](#) [1 图像](#)

概述

产品名称	人SAA ELISA试剂盒(SAA1)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 0.5 ng/ml
范围	0.41 ng/ml - 300 ng/ml
回收率	100 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	93.75	86% - 102%
Serum	114.8	105% - 124%
Plasma	112.1	83% - 137%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Human SAA ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of human SAA in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human Serum Amyloid A coated on a 96-well plate. Standards and samples are pipetted into the wells and Serum Amyloid A present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Serum Amyloid A antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Serum Amyloid A bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

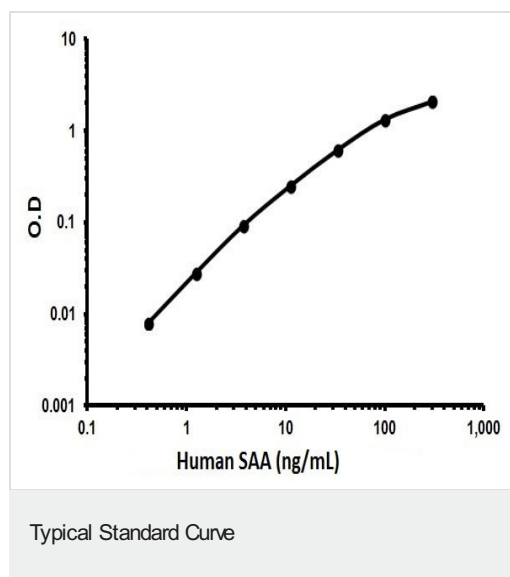
说明 Optimization may be required with urine samples. ab100635 cross reacts with SAA2.
平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
800X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent C	1 x 30ml
Biotinylated anti-Human SAA	2 vials
Recombinant Human SAA Standard (lyophilized)	2 vials
Serum Amyloid A Microplate (12 x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

图片



Representative standard curve using ab100635.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors